

Purity Specification:	≥98%			
Molecular formula:	C19H29N5O3 • XHCl			
CAS number:	Formula weight:	375.47	Expiration date:	22FEB2029

Tests	Results
HPLC	Purity: 100.0 %
IR	Conforms
Mass Spec	[M+H] ⁺ : 376.4
Melting point	239 - 240 °C
TLC	Purity: 100 %
NMR	Conforms
Counterion	Conforms

Reviewed and approved by: Jennifer LaBrecque

Gen Labucque

Copyright Cayman Chemical Company, 10/12/2018

WWW.CAYMANCHEM.COM